

## Anti-CDK4

Rabbit Polyclonal Antibody

Catalog # C31-63R

Lot # B3216-32

### Cited Applications

ELISA, IP, WB

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the 24kDa CDK4 (p34) protein.

### Cross Reactivity

Human, Rat and Mouse

### Host/Isotype/Clone#

Rabbit, Antiserum

### Immunogen

The antibody was produced against synthesized peptide corresponding to the C-Terminus of the human CDK4 (p34) protein conjugated to KLH.

### Formulation

0.01% (w/v) Sodium Azide

### Stability

1yr at -20°C from date of shipment

### Scientific Background

CDK4 is a member of the cyclin-dependent protein kinase family and is involved in the control of cell proliferation during the G1 phase of cell cycle. CDK4 forms a complex with the D-type cyclins and is inhibited by p16 (cyclin-dependent kinase inhibitor-2). CDK4 can mediate phosphorylation of the C-terminal region of RB protein leading to an active transcriptional repression of E2F complex (1). CDC37 and HSP90 can preferentially associate with the fraction of CDK4 not bound to D-type cyclins. SMAD3 is a major physiologic substrate of the G1 cyclin-dependent kinases CDK4 and CDK2 (2).

### References:

1. Harbour, J W. et al: Cdk phosphorylation triggers sequential intramolecular interactions that progressively block Rb functions as cells move through G1. Cell 98: 859-869, 1999.
2. Matsuura, I. et al: Cyclin-dependent kinases regulate the antiproliferative function of Smads. Nature 430: 226-231, 2004.

## Anti-CDK4

Rabbit Polyclonal Antibody

Catalog #	C31-63R
Lot #	B3216-32
Purification	Delipidation and Defibrination
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**